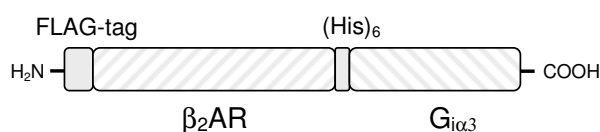


β_2 -AR- $G_{i\alpha 3}$

β_2 -Adrenergic Receptor $G_{i\alpha 3}$ fusion protein
 human, recombinant, Sf9 insect cells

Cat. No.	Amount
PR-541	1 ml



For *in vitro* use only
 Quality guaranteed for 12 months
 Store at -80°C

Avoid freeze / thaw cycles

Form

Membrane suspension. Supplied in 75 mM Tris-HCl
 pH 7.4, 12.5 mM MgCl₂ and 1 mM EDTA.

Molecular Weight

90 kDa

Activity

7.6 pmol/mg

Description

β_2 -Adrenergic receptor- $G_{i\alpha 3}$ is a fusion protein in which the $G_{i\alpha 3}$ N-terminus is linked to the β_2 -adrenoceptor (β_2 AR) C-terminus via a hexahistidine (His₆)-tag.

The β_2 AR is activated by the catecholamine epinephrine and couples to the G-protein G_s to mediate adenylate cyclase (AC) activation. β_2 ARs are found in numerous tissues and cell types including vascular and bronchial smooth muscle cells, leukocytes and liver. β_2 ARs mediate smooth muscle relaxation, inhibition of leukocyte function and activation of glycogenolysis.

In addition to G_s -proteins, the β_2 AR can also couple to G_T -proteins. The β_2 AR- $G_{i\alpha 3}$ fusion protein ensures a defined 1:1 stoichiometry of the receptor and the $G_{i\alpha 3}$ subunit. Coupling efficiency in the β_2 AR- $G_{i\alpha 3}$ fusion protein is lower than in β_2 AR- $G_{s\alpha}$ fusion proteins (cat.# PR-532 and PR-544) as assessed by high-affinity agonist binding and [³⁵S]GTP γ S binding.

The β_2 AR contains a N-terminal FLAG-tag® for immunochemical detection.

Selected References:

- Asano *et al.* (1984) Activation of the inhibitory GTP-binding protein of adenylate cyclase, G_i , by β -adrenergic receptors in reconstituted phospholipid vesicles. *J. Biol.Chem* **259**:9351.
- Seifert *et al.* (1998) Reconstitution of β_2 -adrenoceptor-GTP-bindingprotein interaction in Sf9 cells: High coupling efficiency in β_2 -adrenoceptor- $G_{s\alpha}$ fusion protein. *Eur. J. Biochem.* **255**:369.
- Wenzel-Seifert *et al.* (2000) Molecular analysis of β_2 -adrenoceptor coupling to G_s , G_T , and G_q -proteins. *Mol. Pharmacol.* **58**:954.